

Actinomycin D

Intercalating transcription inhibitor

An intercalating transcription inhibitor and antineoplastic antibiotic. Inhibits transcription by binding to DNA at dG residues. Induces apoptosis in many cell lines, but suppresses in PC12 cells induced by topoisomerase II inhibitor etoposide. Used at 5-10 µg/ml in cell culture for halting transcription. The product is not sterile.

Citations: 20

[View Online »](#)

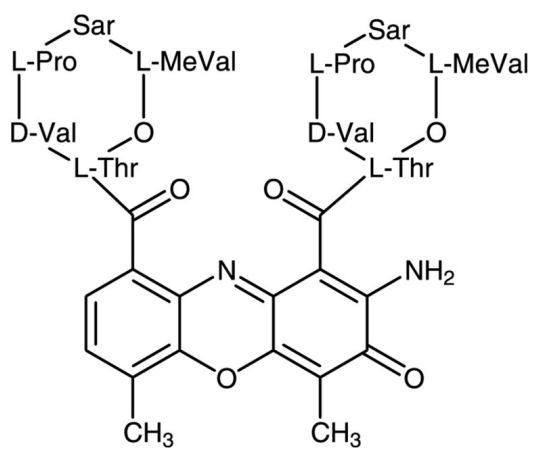
Ordering Information

[Order Online »](#)

BML-GR300-0005	5mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Sensitive to strong acids and bases. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Handling	Protect from light. Hygroscopic.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Red-orange crystalline solid.
CAS	50-76-0
Formula	$C_{62}H_{86}N_{12}O_{16}$
MW	1255.5
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (>20mg/ml), 100% ethanol (1mg/ml) or methanol.
Source	Isolated from <i>Streptomyces parvulus</i> .

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)